

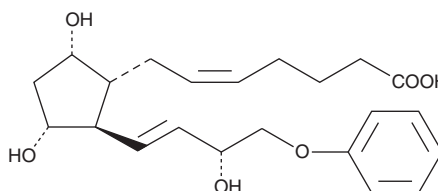
# PRODUCT INFORMATION



## 16-phenoxy tetranor Prostaglandin F<sub>2α</sub>

Item No. 16760

**CAS Registry No.:** 51705-19-2  
**Formal Name:** (5Z)-7-[(1R,2R,3R,5S)-3,5-dihydroxy-2-[(1E,3R)-3-hydroxy-4-phenoxy-1-buten-1-yl]cyclopentyl]-5-heptenoic acid  
**Synonym:** 16-phenoxy tetranor PGF<sub>2α</sub>  
**MF:** C<sub>22</sub>H<sub>30</sub>O<sub>6</sub>  
**FW:** 390.5  
**Purity:** ≥98%  
**UV/Vis.:** λ<sub>max</sub>: 219 nm  
**Supplied as:** A 10 mg/ml solution in methyl acetate  
**Storage:** -20°C  
**Stability:** ≥4 years



Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

### Laboratory Procedures

16-phenoxy tetranor Prostaglandin F<sub>2α</sub> (16-phenoxy tetranor PGF<sub>2α</sub>) is supplied as a solution in methyl acetate. To change the solvent, simply evaporate the methyl acetate under a gentle stream of nitrogen and immediately add the solvent of choice. Solvents such as ethanol, DMSO, and dimethyl formamide purged with an inert gas can be used. The solubility of 16-phenoxy tetranor PGF<sub>2α</sub> in these solvents is approximately 100 mg/ml.

16-phenoxy tetranor PGF<sub>2α</sub> is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, the methyl acetate solution of 16-phenoxy tetranor PGF<sub>2α</sub> should be diluted with the aqueous buffer of choice. The solubility of 16-phenoxy tetranor PGF<sub>2α</sub> in PBS (pH 7.2) is approximately 0.8 mg/ml. We do not recommend storing the aqueous solution for more than one day.

### Description

16-phenoxy tetranor PGF<sub>2α</sub> is a metabolically stable analog of PGF<sub>2α</sub>. It binds to the FP receptor on ovine luteal cells with much greater affinity (440%) than PGF<sub>2α</sub>.<sup>1</sup>

### Reference

1. Balapure, A.K., Rexroad, C.E., Jr., Kawada, K., *et al.* Structural requirements for prostaglandin analog interaction with the ovine corpus luteum prostaglandin F<sub>2α</sub> receptor. *Biochem. Pharmacol.* **38(14)**, 2375-2381 (1989).

#### WARNING

THIS PRODUCT IS FOR RESEARCH ONLY - NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

#### SAFETY DATA

This material should be considered hazardous until further information becomes available. Do not ingest, inhale, get in eyes, on skin, or on clothing. Wash thoroughly after handling. Before use, the user must review the [complete](#) Safety Data Sheet, which has been sent via email to your institution.

#### WARRANTY AND LIMITATION OF REMEDY

Buyer agrees to purchase the material subject to Cayman's Terms and Conditions. Complete Terms and Conditions including Warranty and Limitation of Liability information can be found on our website.

Copyright Cayman Chemical Company, 04/21/2026

#### CAYMAN CHEMICAL

1180 EAST ELLSWORTH RD  
ANN ARBOR, MI 48108 · USA

**PHONE:** [800] 364-9897  
[734] 971-3335

**FAX:** [734] 971-3640

CUSTSERV@CAYMANCHEM.COM  
WWW.CAYMANCHEM.COM